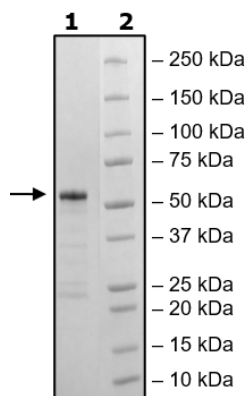


Product Information

Description:	Recombinant human FGFR4 (fibroblast growth factor receptor 4) encompassing amino acids 460-802(end). This construct contains an N-terminal GST-tag. The recombinant protein was affinity purified.
Species	Human
Construct:	FGFR4 (GST-460-802(end))
Concentration:	0.27 mg/ml
Expression System:	Sf9
Purity:	80%
Format:	Aqueous buffer solution.
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 20% glycerol, 3 mM DTT, and 8 mM glutathione
MW:	65 kDa
Genbank Accession:	NM_002011
Stability:	At least 6 months at -80°C.
Storage:	-80°C
Instructions for Use:	Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.
Specific Activity:	50 pmol/min/μg
Assay Conditions:	Activity was measured using a Kinase Buffer (50 mM Tris, pH 7.4, 20 mM MgCl ₂ , 0.1 mg/ml BSA, and 1 mM DTT) containing 20 μM ATP and 0.2 mg/ml poly-(Glu-Tyr) for 45 min at 30°C. Enzyme activity was measured using Kinase-Glo™ plus reagents (Promega).
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



FGFR4 Activity

